

## Anti-ERK2

Rabbit Polyclonal Antibody

**Catalog # M28-63R**

Lot # L020-4

### Cited Applications

- Western blot (1:2500)

*Ideal working dilutions for each application should be empirically determined by the investigator.*

### Specificity

Recognizes only the ERK2 protein

### Cross Reactivity

- ERK2 from human, rat, bovine, mouse, and rabbit cells

*ERK2 from other species may also be detectable*

### Host

Rabbit

### Immunogen

Human ERK2 c-terminal peptide-KLH conjugate

### Formulation

0.1M TRIS pH 8.0, 0.1M glycine

### Stability

Store at 4°C (add 0.1% NaN<sub>3</sub>) for several months, and at -20°C for longer periods. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For optimal performance, avoid repeated handling and multiple freeze/thaw cycles.

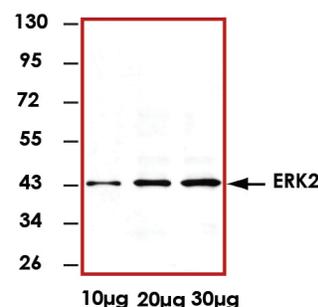
### Scientific Background

ERK2 is a protein serine/threonine kinase that is a member of the extracellular signal-regulated kinases (ERKs) which are activated in response to numerous growth factors and cytokines (1). Activation of ERK2 requires both tyrosine and threonine phosphorylation that is mediated by MEK. ERK2 is ubiquitously distributed in tissues with the highest expression in heart, brain and spinal cord. Activated ERK2 translocates into the nucleus where it phosphorylates various transcription factors (e.g., Elk-1, c-Myc, c-Jun, c-Fos, and C/EBP beta).

### References

1. Boulton, TG. et al: Purification and properties of extracellular signal-regulated kinase 1, an insulin-stimulated microtubule-associated protein 2 kinase. *Biochemistry*. 1991 Jan 8;30(1):278-86.

### Sample Data



Representative western blot with Anti-ERK2 (1:2500) using 5 µg, 10 µg, and 30 µg of HEK-293 cell lysate.

## Anti-ERK2

Rabbit Polyclonal Antibody

Catalog Number

M28-63R

Specific Lot Number

L020-4

Purification

Affinity purified using Protein G

Concentration

0.25 µg/µL

Stability

1 yr At -20°C from date of shipment

Storage & Shipping

Store product at -20°C. For optimal storage, aliquot antibody into smaller quantities after centrifugation and store at recommended temperature. For optimal performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on ice packs.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: [orders@signalchem.com](mailto:orders@signalchem.com)  
[www.signalchem.com](http://www.signalchem.com)

**FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.**